

# PRODUCT INFORMATION

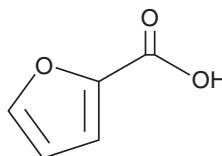


## 2-Furoic Acid

Item No. 45449

**CAS Registry No.:** 88-14-2  
**Formal Name:** 2-furancarboxylic acid  
**Synonyms:** 2-Carboxyfuran, Furan-2-carboxylic Acid, NSC 8842, NSC 58965, Pyromucic Acid

**MF:** C<sub>5</sub>H<sub>4</sub>O<sub>3</sub>  
**FW:** 112.1  
**Purity:** ≥98%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

2-Furoic acid is supplied as a solid. A stock solution may be made by dissolving the 2-furoic acid in the solvent of choice, which should be purged with an inert gas. 2-Furoic acid is soluble (≥10 mg/ml) in DMSO and ethanol. 2-Furoic acid is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2).

### Description

2-Furoic acid is a synthetic and biosynthetic intermediate.<sup>1,2,3</sup> It is a synthetic intermediate in the conversion of lignocellulose-derived furural into 2,5-furandicarboxylic acid, a compound used in the synthesis of pharmaceuticals, pesticides, and polymers.<sup>1</sup> 2-Furoic acid has been used as an organic ligand in metal-organic coordination complexes and coordination polymers.<sup>2</sup> It is also an intermediate in the furfural degradation pathway in bacteria.<sup>3</sup>

### References

1. Zhang, Y., Cui, H. and Xia, H. Recent advances in the synthesis of 2-furoic acid and 2,5-furandicarboxylic acid from furfural. *ChemSusChem* **18**(3), e202401390 (2025).
2. Ejarque, D., Sánchez-Férez, F., Félez-Guerrero, N., *et al.* Pyridine-driven assembly of Zn(II) and Cd(II) complexes with 2-furoic acid. The role of water in a structural transformation. *CrystEngComm* **25**(18), 2739-2754 (2023).
3. Wierckx, N., Koopman, F., Ruijsenaars, H. J., *et al.* Microbial degradation of furanic compounds: biochemistry, genetics, and impact. *Appl. Microbiol. Biotechnol.* **92**(6), 1095-1105 (2011).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/13/2026

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM